

Alexa Fluor® 647 Microscale Protein Labeling Kit (A30009)

Quick Facts

Storage upon receipt:

Component A

- 4°C or ≤–20°C, if desired
- Protect from light
- Protect from moisture

Component B

- Protect from moisture

Components C and D

- Store at room temperature

Component E

- 2–6°C only; DO NOT FREEZE

Abs/Em of conjugate: 650/668 nm

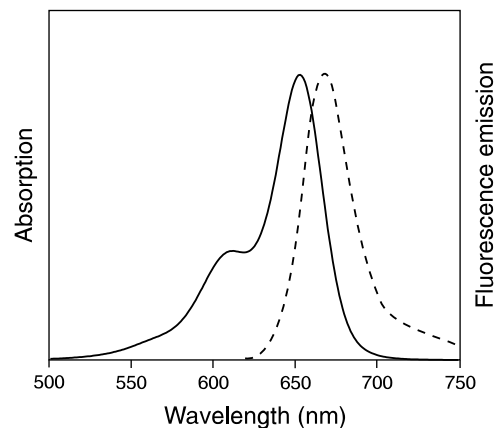


Figure 1. Absorption (—) and fluorescence emission (---) spectra of Alexa Fluor 647 goat anti-mouse IgG antibody in pH 7.2 buffer.

proteins have excitation and fluorescence emission maxima of approximately 650 and 668 nm, respectively (Figure 1). The wide spectral separation of this dye from commonly used red fluorophores such as the Texas Red® and Alexa Fluor 594 dyes makes it an ideal choice for multicolor applications. Moreover, the fluorescence of the Alexa Fluor 647 dye is independent of pH between pH 4 and 10, and conjugates labeled with this dye are typically brighter than similar proteins labeled with Cy®5 dye. Unlike Cy5 dye, the Alexa Fluor 647 dye is resistant to quenching at high degrees of labeling.¹ Each of the vials of Alexa Fluor 647 succinimidyl ester supplied in the kit is sufficient for labeling one sample (20–100 µg at a concentration of 1 mg/ml) of protein that has a molecular weight between ~12 and ~150 kDa.

Introduction

The Alexa Fluor® 647 Microscale Protein Labeling Kit provides a convenient means for labeling small amounts (20–100 µg) of purified protein with our superior Alexa Fluor 647 dye. This kit has been optimized for labeling proteins with molecular weights between 12 and 150 kDa, and contains everything needed to perform three labeling reactions and to separate the resulting conjugates from excess dye. Convenient spin columns are used to purify the labeled protein with yields between 60 and 90%, depending primarily on the molecular weight of the starting material. Labeling and purification can be completed in as little as 30 minutes. For labeling larger amounts of protein, we recommend either the easy-to-use Alexa Fluor 647 Protein Labeling Kit (A20173), which is optimized for 1 mg samples of >40 kDa proteins, or the Alexa Fluor 647 Monoclonal Antibody Labeling Kit (A20186), which is optimized for 100 µg samples of mono- or polyclonal antibodies.

The Alexa Fluor 647 reactive dye has a succinimidyl ester moiety that reacts efficiently with primary amines of proteins to form stable dye-protein conjugates. Alexa Fluor 647 dye-labeled

Materials

Kit Contents

- **Alexa Fluor 647 carboxylic acid, succinimidyl ester** (Component A) 3 vials
- **Sodium bicarbonate** (Component B) 84 mg
- **Reaction tubes** (Component C) 3
- **Spin columns** (Component D) Nanosep® MF 0.2 µm centrifugation devices, 3
- **Purification resin** (Component E) Bio-Gel® P-6 fine resin suspended in PBS,* ~3 ml settled

Storage and Handling

Upon receipt, all kit components can be stored refrigerated at 2–6°C until required for use. The reactive dye (Component A) may be stored at ≤–20°C, if desired, and should be protected from light. Components A and B should be protected from moisture. DO NOT FREEZE COMPONENT E. When stored appropriately, the kit components should be stable for approximately 6 months.

Protein Preparation

IMPORTANT: The purified protein should be at a concentration of 1 mg/ml in a buffer that does not contain primary amines (e.g., ammonium ions, Tris, glycine, ethanolamine, triethylamine, glutathione), or imidazole. All of these substances will significantly inhibit protein labeling. Also, partially purified protein samples, or protein samples containing carriers like BSA (e.g., antibodies) will not be labeled well and should not be used. The presence of low concentrations (<0.1% (w/v)) of biocides, including sodium azide and thimerosal, will not significantly affect the labeling reaction.

To aid in removing low molecular weight components from the protein sample (desalting) prior to labeling, it is possible to use dialysis or small-scale gel filtration. There are a number of easy-to-use, low-volume dialysis options available, including Tube-O-DIALYZER micro-dialysis cartridges from Genotech (www.gbiosciences.com).

We suggest PBS, pH 7.2–7.5, as a suitable prelabeling dialysis buffer, although 100 mM sodium bicarbonate buffer can also be used. If bicarbonate buffer is used, you may omit step 1.1 of the labeling reaction as well as the addition of 1/10 volume of bicarbonate in step 1.2.

Labeling Reaction

Table 1 shows the recommended amount in nanomoles of reactive dye that should be added for each nanomole of protein to be labeled. This is the dye: protein molar ratio (MR). The MR values are based on two parameters: 1) the molecular weight of seven representative proteins ranging from 12 to 150 kDa, and 2) the optimal Alexa Fluor 647 degree of labeling (DOL) for these proteins, as determined in our laboratories. Because your proteins may differ from those listed, the recommended MR for both a lower and a higher DOL are also included in Table 1. For your initial labeling attempt, choose the optimal MR for the protein listed in Table 1 that is closest in molecular weight to the one you are labeling. Use the lower and higher MR as a guide for relabeling if your protein is under- or overlabeled (see Notes). Table 1 also shows typical % yields for the indicated Alexa Fluor 647 conjugates.

Table 1. Recommended Alexa Fluor 647 dye: protein molar ratios (MR) and typical yields for labeling 12–150 kDa proteins.

Protein (MW in kDa)	For Lower DOL	For Optimal DOL	For Higher DOL	% Yield
parvalbumin (12)	≤6	7	≥9	60
soybean trypsin inhibitor (20)	≤6	10	≥12	60–80
thrombin (40)	≤6	12	≥18	70
streptavidin (53)	≤6	12	≥15	84–92
transferrin (80)	≤8	12	≥17	69–81
F(ab) ₂ (100)	≤9	15	≥16	78–84
IgG (150)	≤10	20	≥28	79–97

Use equation 1 to calculate the appropriate volume of reactive dye stock solution to use:

Equation 1

$$\frac{[(\mu\text{g protein/protein MW}) \times 1,000] \times \text{MR}}{7.94} = \mu\text{l reactive dye to add to sample}$$

where $\mu\text{g protein}$ is the mass of protein you want to label, protein MW is the molecular weight of your protein in Da, MR is the dye: protein molar ratio from Table 1, and 7.94 is the concentration of the reactive dye stock solution (see step 1.3 below). For example, to label 60 μg of IgG (MW 150,000):

$$\frac{[(60/150,000) \times 1,000] \times 20}{7.94} = 1.0 \mu\text{l of dye}$$

Do NOT prepare the reactive dye stock solution (step 1.3) until you are ready to start the labeling reactions. This reactive dye hydrolyzes in water and therefore should be used immediately.

1.1 Prepare a 1 M sodium bicarbonate solution by adding 1 ml deionized water to the vial of sodium bicarbonate (Component B). Vortex or pipet up and down until the reagent is fully dissolved. The bicarbonate solution will have a pH of ~8.3 and can be stored at 4°C for up to two weeks. It can also be frozen for long-term storage.

1.2 Transfer 20–100 μl of a 1 mg/ml solution of protein (20–100 μg) to a reaction tube (Component C). Add 1/10 volume (2–10 μl) of 1 M sodium bicarbonate, and mix by pipetting up and down several times.

1.3 Add 10 μl dH₂O to one vial of Alexa Fluor 647 succinimidyl ester (Component A). Completely dissolve the contents of the vial by pipetting up and down. The concentration of this reactive dye stock solution is 7.94 nmol/ μl . As noted above, this solution should be prepared immediately before use, and any leftover solution should be discarded.

1.4 Add the appropriate volume of reactive dye solution, based on equation 1, to the reaction tube containing the pH-adjusted protein and mix thoroughly by pipetting up and down several times.

1.5 Incubate the reaction mixture for 15 minutes at room temperature.

Conjugate Purification

2.1 To prepare for separating the labeled protein from unreacted dye, take the container of gel resin (Component E) and one spin column (Component D; Figure 2) out of the kit. Fully resuspend the gel resin by gently rocking the container; do not vortex or use a magnetic stir bar to agitate the material. Fill the upper chamber of the spin column up to the lip with the suspended gel resin; approximately 800 μl of resin will be needed. Centrifuge the spin column at $16,000 \times g$ in a microcentrifuge for a total of 15 seconds (including run-up time). Using a fixed-angle rotor will cause the resin to pack down with a low side and a high side. The edge of the resin bed on the low side should be about 2–3 mm above the green collar, and the edge of the high side should not be above the top lip of the spin column. If a swinging bucket rotor is used, the resin bed should fill about half of the upper chamber (~ 5 mm). Figure 2 illustrates the two parts of the spin column and what the filled spin column should look like after centrifugation. If the bed is too small, add more suspended resin and centrifuge again. If there is too much resin, resuspend the resin in the upper chamber in buffer, remove the necessary amount, and centrifuge again to repack the bed.

2.2 Occasionally, some resin will get into the collection tube during column preparation. When the resin bed is at the correct level, rinse out the collection tube under the spin column several times with buffer to remove any resin particles that may be found there. Replace the resin-containing insert.

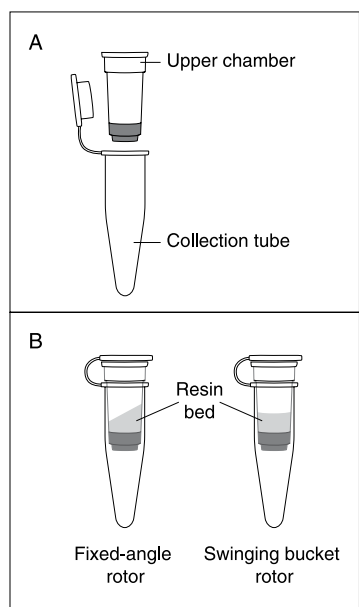


Figure 2. Spin column. (A) An empty column showing the separate parts. (B) Appearance of the resin bed after centrifugation in a fixed-angle rotor or swinging bucket rotor.

If you wish to purify the conjugate in a buffer other than the PBS, pH 7.2, in which the resin is suspended, there are two ways to exchange the buffer. While the resin in the bottle is completely settled, you can decant or aspirate the buffer provided and replace it with another buffer of your choice. Add your buffer to the bottle, mix gently to resuspend the resin, and let it settle completely. Carefully remove the buffer again, and repeat this washing process several times. You can also exchange the buffer after the resin bed is prepared in the spin column, by washing your chosen buffer through the bed several times by brief low-speed centrifugation. The Bio-Gel P-6 fine resin provided is stable between pH 2 and 10.

2.3 After the spin column is prepared, pipet no more than 50 μl of the conjugate reaction mixture onto the center of the resin bed surface. If the volume of conjugate reaction is 51–100 μl , divide it into two aliquots and purify them on separate spin columns. Place the spin column(s) in the microcentrifuge with the high side of the resin bed on the outside. Centrifuge at $16,000 \times g$ for a total of 1 minute.

Note: We have not carefully evaluated reusing a spin column for two aliquots of a larger (≥ 50 μl) sample, and we do not recommend it.

2.4 Each collection tube now contains purified dye-labeled protein in approximately 60–100 μl of buffer. The unreacted dye is retained in the column and the resin will have a blue color (see Notes). The procedure described in steps 1.1–2.3 can be performed in as little as 30 minutes.

Determination of Degree of Labeling (DOL)

Several spectrophotometric methods are available for determining the DOL of the Alexa Fluor 647-labeled protein conjugate. They are based on obtaining the protein concentration by absorbance at 280 nm (A_{280}) and at 650 nm (A_{650}).

3.1 The easiest way to analyze the conjugates spectrophotometrically is using a NanoDrop ND-1000 spectrophotometer (NanoDrop Technologies, Rockland, Delaware, USA). This instrument requires only 2 μl of sample and is a full-featured UV/Vis instrument. A variety of cuvettes are available for use with small sample volumes if you would prefer to not dilute your labeled protein in order to use standard 0.5 or 1.5 ml cuvettes. Quartz cuvettes from Starna Cells, Inc. (Atascadero, California, USA) that hold 15–115 μl of sample are ideal for this purpose.

3.2 The conjugate samples (appropriately diluted) can also be placed in wells of a microplate and read at 280 and 650 nm in a microplate reader that permits the user to specify the desired detection wavelengths, such as the Tecan Safire (Tecan US, Research Triangle Park, North Carolina, USA).

3.3 The conjugate samples can be diluted as necessary prior to measurement of A_{280} and A_{650} using cuvettes and spectrophotometers of your choice. However, excessive dilution of some proteins with low intrinsic A_{280} may prevent you from deriving accurate A_{280} values for your samples. The entire conjugate sample, or a

portion of it, should be diluted only to the minimum volume necessary for your cuvettes and spectrophotometer to avoid relying on readings below the optimal range for your instrument.

3.4 Calculate the concentration of the protein in the sample:

Equation 2

$$\text{Protein concentration (mg/ml)} = \frac{[A_{280} - 0.03(A_{650})] \times \text{dilution factor}}{A_{280} \text{ of protein at 1 mg/ml}}$$

In this equation, 0.03 is a correction factor for the fluorophore's contribution to the A_{280} .

Equation 3

$$\text{Protein concentration (M)} = \frac{\text{answer from Eq. 2}}{\text{protein molecular weight (Da)}}$$

If you know the molar extinction coefficient (ϵ , in $\text{cm}^{-1}\text{M}^{-1}$) of your protein at 280 nm, use this value as the divisor in equation 2 to directly calculate the molarity of the protein.

Example:

You have labeled IgG (MW 150,000; A_{280} at 1 mg/ml = 1.4; ϵ = 210,000 $\text{cm}^{-1}\text{M}^{-1}$) and diluted the sample 1:10 to make the measurements. The readings obtained are $A_{280} = 0.15$ and $A_{650} = 0.90$.

$$\text{Protein concentration (mg/ml)} = \frac{[0.15 - 0.03(0.90)] \times 10}{1.4} = 0.88 \text{ mg/ml}$$

$$\text{Protein concentration (M)} = \frac{0.88}{150,000} = 5.9 \times 10^{-6} \text{ M}$$

$$\text{OR } \frac{[0.15 - 0.03(0.90)] \times 10}{210,000} = 5.9 \times 10^{-6} \text{ M}$$

3.5 Calculate the degree of labeling:

Equation 4

$$\text{DOL} = (\text{moles dye})/(\text{mole protein}) = \frac{A_{650} \times \text{dilution factor}}{239,000 \times \text{protein concentration (M)}}$$

where 239,000 $\text{cm}^{-1}\text{M}^{-1}$ is the approximate molar extinction coefficient of the Alexa Fluor 647 dye.

Example:

$$\text{DOL} = \frac{0.90 \times 10}{239,000 \times 5.9 \times 10^{-6} \text{ M}} = 6.4$$

Note that low molecular weight Alexa Fluor 647-labeled proteins with DOL values of 1–2 may appear very pale blue in color. This does not necessarily mean that the protein did not label well. Alexa Fluor 647 is a very bright dye with a high molar extinction coefficient, and the color of the conjugate may not be a good indicator of conjugation success.

3.6 If the A_{280} of your protein is too low to measure accurately, you can get an approximate protein concentration by estimating the % yield of the conjugate. The % yield from the spin columns is related to the molecular weight of the protein, and typical yields are provided in Table 1. Once you have a % yield,

the approximate protein concentration can be calculated using equation 5:

Equation 5

$$\text{Concentration (mg/ml) labeled protein} = \frac{(\text{mass (mg) starting protein} \times \% \text{ yield})}{\text{volume (ml) recovered}}$$

3.7 Divide the concentration (in mg/ml) by the protein's molecular weight (in Da) to calculate the approximate molar concentration of the protein. To determine the approximate DOL, you still must spectrophotometrically determine the A_{650} of the conjugate. However, this value is usually significantly higher than the A_{280} even if the conjugate has been diluted significantly. The A_{650} is used in equation 4 (step 3.5) to determine the DOL.

Storage of Conjugates

We typically store labeled proteins at 2–6°C, protected from light. It may be necessary to add a stabilizer like BSA (1–10 mg/ml) or glycerol to your conjugate to improve stability. In the presence of 2 mM sodium azide or other biocides, a typical antibody conjugate should be stable at 2–6°C for several months. Your proteins may have special storage requirements. If it is appropriate to do so with your proteins, you can divide the conjugate into small aliquots and freeze them at $\leq -20^\circ\text{C}$ for longer storage. AVOID REPEATED FREEZING AND THAWING, AND PROTECT FROM LIGHT.

Notes

Many protein- and dye-specific properties determine how efficiently a protein can be labeled with an amine-reactive dye. Important factors include the number of solvent-accessible primary amines in the protein, the protein's pI, and its solubility and stability at pH 8–8.3. Reactive labels vary in amine reactivity, often in a protein-specific way, and their behavior can be predicted with confidence for only a few proteins such as antibodies and streptavidin. Thus, the recommended dye: protein molar ratios in Table 1 are given as guidelines only, and we cannot guarantee that they will yield optimal labeling with your particular protein(s).

The number of reactive fluorescent dyes that can be attached to a protein before fluorescence quenching or protein inactivation or precipitation occurs is roughly proportional to the protein's molecular weight. For example, the optimal DOL would usually be 1–2 for a ~20 kDa protein, while the optimal DOL for a ~150 kDa protein, e.g., an IgG, would usually be 5–8. The DOL that you obtain with a protein using the Alexa Fluor 647 Microscale Protein Labeling Kit may be higher or lower than the generally accepted optima. We highly recommend that you evaluate your protein conjugate in its intended application before you conclude that it is under- or overlabeled. A number of conditions can cause under- or overlabeling.

Underlabeling

- Even trace amounts of primary amine-containing components (e.g., Tris, glycine, ammonium ions, ethanolamine, triethylamine, or glutathione) or imidazole in the starting protein sample will

decrease labeling efficiency. The ElutaTube microdialysis vials provided can be used to remove these low molecular weight substances from the protein sample prior to labeling.

- Efficient labeling will probably not occur if the concentration of protein starting material is <1 mg/ml.
- The addition of sodium bicarbonate (step 1.2) is designed to raise the pH of the reaction mixture to ~8, as succinimidyl esters react most efficiently with primary amines at slightly alkaline pH. If the protein solution is strongly buffered at a lower pH, the addition of 1/10 volume of bicarbonate solution will not raise the pH to the optimal level. Either more bicarbonate can be added, or the buffer can be exchanged with PBS, pH 7.2 (and bicarbonate solution added again), or with 100 mM sodium bicarbonate buffer, pH 8.3, by dialysis or another method prior to starting the labeling reaction.
- Because proteins react with fluorophores at different rates and retain biological activity at different degrees of dye labeling, the recommendations shown in Table 1 may not always result in optimal labeling. To increase the DOL, the same protein sample can be relabeled, or a new protein sample can be labeled using more reactive dye. Three vials of reactive dye are provided to allow three labeling reactions. Although this kit was designed for optimal labeling in 15 minutes at room temperature, higher DOL may be obtained with longer incubation times. We have not evaluated incubation times >15 minutes.
- Low molecular weight Alexa Fluor 647-labeled proteins with DOL values of 1–2 may appear very pale blue in color. This does not necessarily mean that the protein did not label well. Alexa Fluor 647 is a very bright dye with a high molar extinction coefficient, and the color of the conjugate may not be a good indicator of conjugation success.
- Underlabeling may be the reason for the fluorescent signal being lower than expected in your application. Should this

* PBS = phosphate-buffered saline (10 mM potassium phosphate, pH 7.2, 150 mM NaCl)

† MWCO = molecular weight cutoff

occur, relabel the sample, or label another sample with more reactive dye.

- We have observed that dye-labeling of some proteins to any degree can destroy their biological activity.

Overlabeling

- Overlabeling may be indicated by the formation of a blue precipitate in the reaction mixture or deposition of blue particles on the upper surface of the resin bed after centrifugation of the conjugate. Precipitation will usually result in a decreased yield of conjugate. If your % yield is <50%, it is likely that the protein is overlabeled. Repeat the labeling reaction with less reactive dye. Some proteins cannot be labeled with amine-reactive dyes under any circumstances and may irreversibly precipitate.
- If no visible precipitate forms during labeling but the fluorescent signal in your application is lower than expected, the fluorescence of the conjugate may be quenched due to overlabeling. To reduce the DOL, use a smaller amount of reactive dye, or try labeling the protein at a concentration of >1 mg/ml. (Note: We have not evaluated labeling efficiency with this kit on proteins at concentrations >1 mg/ml.)
- One cause of apparent overlabeling is inefficient removal of unreacted dye. Although using the spin columns in this kit exactly as described removed all traces of free dye from all of the proteins we tested, it is possible that some free dye may be present in your sample after the purification step. The presence of free dye, which can be determined by thin layer chromatography, will result in erroneously high calculated DOL values. Free dye remaining after use of the spin column can be removed by applying the conjugate to another spin column or by extensive dialysis. Applying no more than 50 µl of conjugate to each spin column is the best way to avoid contamination with free dye.
- We have observed that dye-labeling of some proteins to any degree can destroy their biological activity.

Reference

1. J Histochem Cytochem 51, 1699 (2003).

Product List *Current prices may be obtained from our Web site or from our Customer Service Department.*

Cat #	Product Name	Unit Size
A30009	Alexa Fluor® 647 Microscale Protein Labeling Kit *for 20-100 µg protein* *3 labelings*	1 kit
A20186	Alexa Fluor® 647 Monoclonal Antibody Labeling Kit *5 labelings*	1 kit
A20173	Alexa Fluor® 647 Protein Labeling Kit *3 labelings*	1 kit

Contact Information

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly from Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Please visit our website — probes.invitrogen.com — for the most up-to-date information.

Molecular Probes, Inc.

29851 Willow Creek Road, Eugene, OR 97402
Phone: (541) 465-8300 • Fax: (541) 335-0504

Customer Service: 6:00 am to 4:30 pm (Pacific Time)

Phone: (541) 335-0338 • Fax: (541) 335-0305 • probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209 • Order Fax: (800) 438-0228

Technical Service: 8:00 am to 4:00 pm (Pacific Time)

Phone: (541) 335-0353 • Toll-Free (800) 438-2209
Fax: (541) 335-0238 • probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd.
3 Fountain Drive
Inchinnan Business Park
Paisley PA4 9RF, UK
Phone: +44 (0) 141 814 6100 • Fax: +44 (0) 141 814 6260
Email: euroinfo@invitrogen.com
Technical Services: eurotech@invitrogen.com

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License

For research use only. Not intended for any animal or human therapeutic or diagnostic use. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402. Tel: (541) 465-8300. Fax: (541) 335-0504.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation ® are registered with the U.S. Patent and Trademark Office.

Copyright 2005, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.